

Certificate of Analysis

XP Media CHO Growth A without L-Gln K8860

| | |
|----------------|--------------------|
| Batch # | 99914190901 |
|----------------|--------------------|

| | |
|----------------------------------|---|
| Full Description | XP Media CHO Growth A without L-glutamine |
| Appearance | Light pink solution |
| Components | No material of animal origin is used in the manufacture of this product |
| Volume | 1000 ml of media |
| Sterility | Pass. Note: USP <71> Direct Inoculation method used. |
| Preservatives/Stabilisers | None |
| Endotoxin Levels | <.05 EU/ml endotoxin unit (EU) per ml as determined by LAL method |
| Functional Growth | Supports growth of in house cell line |

This product has been manufactured to ISO9001:2015 Quality Assurance guidelines

For research use only. Store at 2-8°C

Signed: *Kieu Khiem Nguyen*

Date: 28 December 2019

Production Biologist